

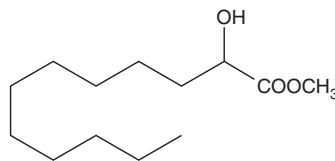
PRODUCT INFORMATION



2-hydroxy Lauric Acid methyl ester

Item No. 24587

CAS Registry No.: 51067-85-7
Formal Name: 2-hydroxy-dodecanoic acid, methyl ester
Synonyms: 2-hydroxy Methyl dodecanoate,
2-hydroxy Methyl laurate,
SFE 13:0;O
MF: C₁₃H₂₆O₃
FW: 230.3
Purity: ≥98%
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

2-hydroxy Lauric acid methyl ester is supplied as a solid. A stock solution may be made by dissolving the 2-hydroxy lauric acid methyl ester in the solvent of choice, which should be purged with an inert gas. 2-hydroxy Lauric acid methyl ester is soluble in organic solvents such as methanol, chloroform, and ethyl ether.

Description

2-hydroxy Lauric acid methyl ester is a hydroxylated fatty acid methyl ester that has been found in coliform bacteria isolated from human sewage.¹

Reference

1. Haznedaroglu, B.Z. and Duran, M. Fatty acid methyl ester (FAME) profiles of total coliforms as a new tool for predicting sources of fecal contamination. *Proc. Water Environ. Fed.* **2005(81-90)**, 7142-7149 (2005).

WARNING
THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA
This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the [complete](#) Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY
Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 03/29/2024

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM